



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Human GIP (active) Calibrator

Catalog Number: C016N-2

Lot Number: A00U0072

Expiration Date: 31 JUL 2023

Storage Condition: 2-8°C

Analyte Description	Amount of Analyte in Calibrator	Units	Stock Concentration of Reconstituted Calibrator (Vol = 250 µL)	Units
hu GIP (active)	7300	pg	29 200	pg/mL

To calculate the **Calibrator Standard 1 (top of curve) concentration**, please refer to the dilution-factor used in the **Prepare Calibrator Standards** section of the respective product insert. Divide the stock concentration of the reconstituted calibrator by this dilution factor to generate the working calibrator standard 1 concentration.

Statement: The above calibrator lot is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	29 AUG 2018